



CERTIFICATE OF ANALYSIS

COA - 11113

Product Description: MSD Blocker A

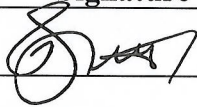

Catalog #/Size: R93BA-2 / 250mL

Lot Number: Y0180905

Expiration Date: 30 NOV 2014

Statement: This product was manufactured and tested for interferents according to approved MSD documents.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		10/1/12
Review/Approval	Michael Eager	Quality Control		14 NOV 2013